

PRODUCT INFORMATION

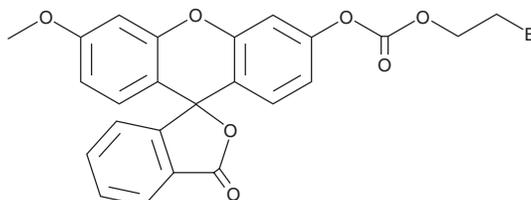


CAY10731

Item No. 26548

CAS Registry No.: 2119597-33-8
Formal Name: carbonic acid, 2-bromoethyl 6'-methoxy-3-oxospiro[isobenzofuran-1(3H),9'-[9H]xanthen]-3'-yl ester

MF: C₂₄H₁₇BrO₇
FW: 497.3
Purity: ≥95%
UV/Vis.: λ_{max}: 223, 274 nm
Ex./Em. Max: 485/535 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

CAY10731 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10731 in the solvent of choice. CAY10731 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of CAY10731 in these solvents is approximately 30 mg/ml.

Description

CAY10731 is a fluorescent probe for hydrogen sulfide (H₂S).¹ Upon reaction with H₂S, fluorescein (FITC) is released and its fluorescence can be used to detect H₂S. FITC displays excitation/emission maxima of 485/535 nm, respectively. CAY10731 is selective for H₂S over glutathione (GSH), cysteine, H₂O₃⁻, SO₃⁻, and H₂O₂ as well as tBuO and OH radicals. It has been used to detect H₂S in rat hippocampal slices and nematodes.

Reference

1. Reja, S.I., Sharma, N., Gupta, M., *et al.* A highly selective fluorescent probe for detection of hydrogen sulfide in living systems: In vitro and in vivo applications. *Chemistry* **23(41)**, 9872-9878 (2017).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/03/2019

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM